

# OXI-TEK TBARS

## assay kit

The sensitivity of measuring thiobarbituric acid reactive substances (TBARS) has made this assay the method of choice for screening and monitoring lipid peroxidation, which is a major indication of oxidative stress. Malondialdehyde (MDA) forms a 1:2 adduct with thiobarbituric acid; the adduct can be measured by fluorometry. Biological specimens contain a mixture of TBARS, including lipid hydroperoxides and aldehydes. In practice, TBARS are expressed in equivalent numbers to MDA.

Citations: 26 [View Online »](#)

Ordering Information [Order Online »](#)

ALX-850-2870	160 tests
--------------	-----------

Manuals, SDS & CofA [View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice – Dangerous Good

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Fluorescent detection

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)